

Invitrogen™ iGlow™ Gel Documentation System

USER GUIDE

Catalog Numbers IGLOW and IGLOW220

Publication Number MAN1001999

Revision A



For Research Use Only. Not for use in diagnostic procedures.

ThermoFisher
SCIENTIFIC



Thermo Fisher Scientific | 3747 N. Meridian Road | Rockford, Illinois 61101 USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: MAN1001999 A (English)

Revision	Date	Description
A	8 May 2026	New document for the iGlow™ Gel Documentation System.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF THE DOCUMENT OR THE PRODUCT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

TRADEMARKS: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2026 Thermo Fisher Scientific Inc. All rights reserved.

Contents

■	CHAPTER 1	Product information	6
		Product description	6
		Contents	6
■	CHAPTER 2	Getting started	8
		Out of box experience (OOBE)	8
		Default accounts	8
		Administrator system setup	9
		Instrument configuration (Administrator only)	9
		Configure date and time	9
		Set password policies	10
		Change the public account password	10
		Network connection	10
		Create an instrument account	11
		Follow the steps to set up your instrument account	11
		Password requirements	11
		Guidelines for strong password	11
		Account management	12
		Reset forgotten password (administrator only)	12
		Change password	12
■	CHAPTER 3	Get Connected	14
		Map instrument profile to a network drive	14
		Create a network drive via settings	14
		Connect iGlow directly to thermal printer	15
■	CHAPTER 4	Imaging	16
		Light sources and sample trays	16
		Image capture workflow	16
		Image preview	17
		Options in image preview screen	17
		Options in contrast versions screen	18
		Options for image adjustment	19

■	CHAPTER 5	Gallery	21
		Manage image files in the gallery	21
		Access the gallery	21
		View and sort images	21
		Search and select images	21
		Monitor storage capacity	22
		Open an image	22
■	CHAPTER 6	Save and export image files	23
		Export options	23
		Additional information	24
■	CHAPTER 7	Band excision	25
		Excise bands from gels requiring UV illumination	25
		Excise bands from gels requiring Blue illumination	26
		Excise bands from gels requiring white illumination	27
■	CHAPTER 8	Instrument maintenance	28
		Administrator settings	28
		Standard users settings	30
		Run the instrument manual self-test	30
		Clean the instrument	31
■	APPENDIX A	Instrument specifications	32
		Instrument dimensions and specifications	32
		Electrical requirements	32
		Environmental requirements	33
		Network requirements	34
		Instrument connections	34

■	APPENDIX B	Safety	35
		Symbols on this instrument	35
		Standard safety symbols	36
		Control and connection symbols	36
		Conformity symbols	36
		Safety information for instruments not manufactured by Thermo Fisher Scientific	38
		Instrument safety	38
		General	38
		Physical injury	39
		Electrical safety	40
		UV safety	41
		Cleaning and decontamination	41
		Instrument component and accessory disposal	41
		Safety and electromagnetic compatibility (EMC) standards	42
		Safety and EMC standards	42
		EMC standards	42
		Environmental design standards	43
		Chemical safety	44
		Biological hazard safety	45
■	APPENDIX C	Documentation and support	46
		Customer and technical support	46
		Limited product warranty	46



Product information

IMPORTANT! Before using this product, read and understand the information in the “Safety” appendix in this document. Wear appropriate protective eyewear, clothing, and gloves.

Product description

The Invitrogen™ iGlow™ Gel Documentation System enables users to image nucleic acid and protein gels stained with fluorescent and visible stains. The iGlow™ system features intuitive workflows and delivers high-resolution images. The built-in touchscreen interface, three distinct LED transilluminator light sources (blue, UV, and white), and scientific CMOS camera enable fast, high-quality data capture. To analyze data, the instrument can be integrated with a license free standalone iBright™ Analysis Software <https://www.thermofisher.com/us/en/home/technical-resources/software-downloads/ibright-western-imager.html>.

Contents

The iGlow™ Gel Documentation System is supplied with the following standard components:

Table 1 iGlow™ Gel Documentation System – Shipment Contents

Category	Components list	Quantity	Catalog number
Instrument	iGlow™ Gel Documentation System (110V or 220V-region specific)	1	IG0001 or IG0002
Accessories ^[1]	UV Tray iGlow™	1	IGA011
	Blue Tray iGlow™	1	IGA012
	White Tray iGlow™	1	IGA013
	Safety Guard (UV SHIELD iGlow™)	1	IGA015
	Tray Holder iGlow™	1	IGA014
	Power cord (region specific)	1	–
Documents	Packaging Checklist	1	–
	Quick Reference Guide	1	–

Table 1 iGlow Gel Documentation System – Shipment Contents *(continued)*

Category	Components list	Quantity	Catalog number
Optional Accessories ^[2]	Thermal Printer	1	A45235
	USB Wi-Fi Module	1	A26774

^[1] The following items, except the power cord, can be purchased separately

^[2] The following items can be purchased separately and are not included in the shipment

This guide helps you set up and start using iGlow™ to streamline your workflow and improve efficiency

Out of box experience (OOBE)

Note: The OOBE feature is available only during the initial use.

1. Remove the iGlow™ Gel Documentation System and accessories from the packaging.
2. Place the instrument on a clean, stable, level laboratory surface.
3. Connect the supplied power cord.
4. Turn on the instrument using the top-right power switch.
5. Allow the system to complete initialization.
6. Review and accept the End User License Agreement (EULA).
7. On the **Welcome** screen, sign in with the provided default credentials and follow the setup prompts.

Default accounts

The system includes **two default accounts** that provides access to the imaging platform:

Table 2 Default user accounts

Account type	Username	Password
Administrator	Administrator	AA
Standard user	Public	aa

First-time setup (administrator):

1. Log-in as **Administrator** using the provided default account credentials.
2. At first login, follow the prompt to **update your password**.
3. Create and confirm the new password.
4. Log in again using your updated credentials.
5. Configure the instrument settings as required for the administrator roles.
6. Set the password policy (applies to all users).

Administrator system setup

1. Log in using the Administrator account.
2. Configure system settings, including:
 - a. Time zone
 - b. Date format

Note: We strongly recommend configuring the password policy during installation in accordance with institutional requirements, as it applies to all user accounts. Changing the policy later may require existing users to reset their passwords.

These settings can be updated at any time as needed.

For additional details, see the administrator privileges section (“Administrator settings” on page 28).

Instrument configuration (Administrator only)

After signing-in with the updated password, the **Administrator** is directed to the **Home** screen

Access **Settings** from the **Home** screen to:

Configure date and time

1. From the Settings menu, go to instrument and select **Time**.
2. Configure the following options as needed.
 - Select the appropriate **Time Zone**.
 - Change **Date**.
 - Change the **Time**.
3. Select **Change** to apply the settings.
4. (Optional) Alternatively, select **Synchronize** to automatically update the date and time using an internet connection.

Set password policies

1. From the Settings menu, select **Permissions**.
2. Select **Password Policy**.
3. Configure the password policy according to institutional requirements, including:
 1. Password composition
 2. Password minimum length
 3. Password validity (in months)
 4. Maximum number of password attempts
 5. Previous password retention

Change the public account password

Note: Public account can be used by all users.

See details of password change in the **Account Management** section.

Record the new password securely and follow your institution's guidelines for password management and sharing. Alternatively, select **Save Password** to enable automatic login to the "**Public**" account

Network connection

This section describes how to connect the iGlow™ instrument to a wired or wireless network.

1. From the **Home** screen, touch **Settings**, touch **Instrument** and touch **Network**.
2. Review the network details displayed at the upper-left corner of the screen
 - The instrument ID
 - Subnet Mask
 - IP address
3. To connect the instrument to a wired network:
 - Connect a live ethernet cable to the ethernet port on the rear panel of the instrument
 - Network status icons are displayed in the upper-right corner of the Instrument Settings screen under the Network

Note: DHCP-enabled networks connect automatically.

Note: Contact your local IT department if assistance is required.

4. To connect the instrument to a wireless network:
 - Insert the High-Power USB Wi-Fi Module (Cat. No. [A26774](#)) into an available USB port
 - Wait for the list of available wireless networks to appear
 - Select the desired wireless network
 - Follow the on-screen instructions to complete the connection

Create an instrument account

Users can create individual accounts from the welcome screen. Passwords must meet administrator-defined policies.

Follow the steps to set up your instrument account

1. From the welcome page, select **Create**.
2. Enter a **Username**.
3. Enter a **Password** then re-enter the password to confirm.
4. Select **Create** to complete account creation.
5. When the account is successfully created, the system displays the confirmation message **Account Added**.
6. Return to the **Welcome** page and sign in using the newly created account credentials.
7. (Optional): Enable **Save Password** to store the password on the instrument

Note: If **Save password** is enabled, the account can be accessed by any user of this instrument.

Password requirements

The instrument does not allow commonly used passwords and simple sequences.

If a restricted password is entered, the system displays one of the following messages:

- **Password requirement not met. Press to view** – displays the password criteria
- **Please choose a stronger password**

Guidelines for strong password

A strong password must meet the following criteria based on the configured settings:

- Use a mix of uppercase and lowercase letters
- Include at least one number and one special character
- Avoid commonly used words or phrases, such as password or your username

- Avoid simple or repeated patterns, including:
 - Repeated numbers (for example, 111111)
 - Sequential numbers (for example, 12345678)
 - Adjacent keyboard characters (for example, qwerty)

Account management

Reset forgotten password (administrator only)

1. If a standard user forgets their password, only the administrator can generate a temporary password.
2. The user must create a new password at the next login (see **Change Password**).
3. If the administrator forgets their password, follow the procedure to reset access.
 - a. Contact Technical Support (<https://www.thermofisher.com/store/v2/contact-us/technical-support>) to obtain the electronic key file: SYSTEM LOGIN [instrument serial number] KEY.
 - b. Download the file and copy it to the root directory of a USB drive.
 - c. Insert the USB drive into the instrument.
 - d. On the Welcome page, select **Administrator** and click **Forgot Password?**.
 - e. When prompted, select **Proceed** to initiate the reset.
 - f. On the **Update Password screen**, enter and confirm a new password (per password policy), then select **Update**.
 - g. Confirm successful password update and select **Done**.
 - h. Return to the Welcome page, select **Administrator**, enter the new password (optionally, select **Save Password**), and click **Login**.

Change password

Use this procedure to update your existing password, or to change your password after logging in with a temporary password provided by the administrator.

1. From the Settings menu, select **Permissions**.
2. Select **Account Management**.
3. Select the account to be modified.
4. Select **Edit**.
5. Change the password.

6. Select **Confirm** to apply the changes.

Note: If **Saved Password** is selected at login, any password changes made using this procedure are automatically stored and used for subsequent logins until the password is changed again.



Get Connected

Map instrument profile to a network drive

You can create a Network Drive location in iGlow™ using the **Settings** menu.

IMPORTANT! iGlow™ stores a single set of network (Windows/domain) credentials for all network export locations.

The first user to connect to a network path establishes the shared credentials used for all subsequent network exports.

Other users must request from that first user to add an export locations on their behalf, unless they have access to the same network username and password.

Create a network drive via settings

1. Log in to your user account.
2. Select **Settings**.
3. Navigate to **System**.
4. Select **Export Locations**.
5. Select **Add**.
6. When prompted, enter the **Network Drive path** using back slashes:
 - a. Enter the **IP address**, **server name**, and **file path**.
 - b. Example: \\10.XX.X.XX\shared\R&D_Data\John Smith\2025.
7. Please enter the username and password associated with the network account.
8. Select **Connect** to complete setup.

Connect iGlow directly to thermal printer

The iGlow™ system supports direct USB connection to the recommended thermal printer.

1. Use the recommended Thermal Printer for iGlow Imaging Systems (Cat no [A45235](https://www.thermofisher.com/order/catalog/product/A45235)) (<https://www.thermofisher.com/order/catalog/product/A45235>).
2. Using an A to B type USB cable, connect the instrument (USB-A) directly to the thermal printer (USB-B).
3. Connect the power cord to the printer and plug into a power outlet.
4. Power on the printer.
5. On the iGlow™ instrument, when a captured image is displayed in the viewport, there will be a printer icon located at the bottom of the viewport.
6. Select the icon to print this image on the thermal printer.
7. A pop-up screen will offer options to add the image name to the report and/or to invert the image.

Note: Recommended: UVP High Gloss Thermal Printer Paper (Cat No [UPP89017401](https://www.fishersci.com/shop/products/uvp-thermal-printer-paper/UPP89017401)) (<https://www.fishersci.com/shop/products/uvp-thermal-printer-paper/UPP89017401>).

From the welcome page, select your account and sign-in.

The Home screen displays the available light sources (**UV**, **Blue**, or **White**) provides access to gel excision and to the **Gallery**.

Light sources and sample trays

iGlow™ has three illumination sources that provide flexibility for imaging numerous gel stains. The following table describes how to image commonly available gel stains.

Table 3 Imaging Options by Light Source

Illumination module	Sample tray	Compatible Sample Type
UV	UV Tray	- DNA and RNA gels stained with fluorescent nucleic acid stains (for example, Ethidium Bromide, SYBR™ Safe, SYBR™ Green, SYBR™ Gold and similar products). - Fluorescent-stained protein gels (for example, SYPRO™ Ruby, Coomassie Fluor Orange, Oriole and similar products).
Blue	BLUE tray	- DNA and RNA gels stained with fluorescent nucleic acid stains (for example, Ethidium Bromide, SYBR™ Safe, SYBR™ Green, SYBR™ Gold and similar products). - Fluorescent-stained protein gels (for example, SYPRO™ Ruby, SYPRO™ Red, Coomassie Fluor Orange, Flamingo and similar products).
White	WHITE Tray	Visible-stained protein gels (for example, coomassie blue stains and silver stain).

Image capture workflow

1. Open the drawer.
2. Place the desired sample tray onto the transilluminator box.

3. Use the alignment guides to ensure that the tray is securely and correctly positioned.

Note: The system automatically detects the installed tray and enables the correct illumination source.

4. Place the sample in the center of the tray.
5. Close the drawer.
6. Select the highlighted illumination mode to capture the image.
 - a. The software automatically calculates the optimal exposure time, captures the image, and displays it in the **Image Preview** screen.

IMPORTANT! Do not open the drawer during the image capture process. Opening the drawer will terminate the image capture and return the instrument to the **Home** screen.

Image preview

The **Image Preview** screen allows users to review the captured image and recapture it if needed. Recaptures can be performed using a pre-saved method, manual exposure settings, or by adjusting the field of view (FOV) or light intensity. If no changes are required, the image can be optionally renamed and saved to the selected export location (see **Save and Export Files**).

Options in image preview screen

The following table lists the options available on image preview screen.

Table 4 Image Preview Options

Icons	Description of action
 Rename	Change the file name. Note: Default file naming (<i>YearMonthDayHourMinuteSecondMillisecond</i> (for example, 2025092917330223)).
 Methods	Load predefined acquisition methods. Note: The method information is accessible next to the method name.
 FOV	Select the desired field of view (FOV). Available options are 14 × 21 cm, 10 × 15 cm and 8 × 12 cm Note: The default FOV is defined in the administrator settings.

Table 4 Image Preview Options (continued)

Icons	Description of action
 Light Intensity	Adjust the light intensity to 20%, 40%, 60%, 80%, or 100% for UV or Blue light. Default Settings: - UV: 60% (2–2.5 mW/cm²) - Blue: 60% (13–14 mW/cm²)
 Auto Exposure	Recapture the image with the optimal exposure time.
 Manual Exposure	Recapture the image with a manual set exposure time. Note: The most recent exposure time is displayed.
 Save	Save/export the image to a USB drive or a network drive location.
 Contrast versions	Display the three contrast versions (Low, Medium and High) and a pseudo-colored image.

To adjust the contrast levels of the image, navigate to the **Contrast Versions** screen.

Options in contrast versions screen

From the **Image Preview** screen, select **Contrast Versions** to view the three contrast versions and a pseudo-colored image. This screen provides access to image information, contrast editing tools, file saving options, the image Gallery, and navigation back to the **Home** screen.

Table 5 Contrast Versions sub-options

Icons	Description of action
 Gallery	The on-board Gallery contains all image files captured under the logged-in user profile.
 Edit	Edit a captured image: rotate, flip, straighten, invert, and crop. For details, see Table 6.
 Print	Print the selected contrast version of the image using a thermal printer. Note: iGlow™ is compatible with the Sony UP-X898MD Digital Graphic Printer.

Table 5 Contrast Versions sub-options (continued)

Icons	Description of action
 Export	Generate a PDF report and export it to a USB drive or network drive location.
 Save	Save the image file to a USB drive or a network drive location.
 Histogram	Manually adjust grayscale levels using Histogram slider controls. • Set Range - adjust the intensity range. • Back - restore original range. • Auto Contrast - automatically apply the optimal intensity range. • Saturation - display saturated signals in red.
 Home	Return to the Home screen.
	Image information is displayed at the top corners of each contrast version, including user account details and image acquisition parameters.
	Select a color from the Color bar to apply to the image captured with White illumination.
	Select a color from the Color bar to apply to the image captured with UV and Blue illumination.

Options for image adjustment

Captured images can be further edited for publication purposes.

Note: Modified images can only be saved to external locations (USB drive or network). The original image stored in the **Image Gallery** is not altered.

Table 6 Image Adjustment tools

Icons	Description of action
  Rotate left Rotate right	Rotate the image in 90° increments to the left or right.
  Flip horizontal Flip vertical	Flip the image vertically or horizontally.

Table 6 Image Adjustment tools *(continued)*

Icons	Description of action
 Custom rotate	Rotate the image to a custom angle.
 Invert	Invert the displayed image.
 Crop	Crop the image to remove unwanted areas outside the crop box.



Gallery

Manage image files in the gallery

The on-board **Gallery** contains all image files captured under the logged-in user profile. Image files appear as thumbnails, and users can filter, search, delete, rename, or export files.

Access the gallery

The Gallery can be accessed in two ways:

- From the **Home** screen.
- From the Contrast Version screen, accessed via the **Image Preview** screen.

Note: Each logged-in user has access to their own gallery.

Note: Administrators can access all users galleries.

View and sort images

- The image gallery is organized by acquisition date, with the corresponding date displayed at the top of each section for easy reference.
- Images can be sorted from newest to oldest.
- Use the calendar to adjust the date range for image display.

Note: By default, only images from the last 30 days are displayed. To view older images, adjust the date range using the calendar.

Search and select images

- Images can be searched by name or date.
- To select all images:
 - Select "**Select all**"
 - Select "**Deselect All**" to clear all selections
 - Images can also be selected by date

Monitor storage capacity

- The **total gallery size** and **available disk space** are displayed above the image list.
 - If the available disk space is falls below the preset threshold, the software prompts the user to delete image files before continuing image acquisition.

Note: The disk space threshold that triggers this notification is configured by the **Administrator**.

Open an image

Only one image can be opened at a time.

- Select an image thumbnail
- Double-click the image to open it
- The image opens displaying three contrast versions and a pseudo-colored version



Save and export image files

The Save and export feature allows you to store images within the system or download them for external use. You can export images in supported file formats for reporting, sharing, or further analysis.

Export options

Images can be saved or exported immediately after image capture using one of four available methods. When a save or export option is selected, a pop-up screen is displayed with the following options:

- **File type**
- **Export location**
- Optional actions, including:
 - Renaming the file
 - Exporting image information
 - Saving a **high-resolution** image
 - Saving **all contrast versions** of the image

These options allow users to customize how image files and associated data are stored or exported.

Note: All the files saved will be exported into a folder labeled with the export date (for example, 20260206).

Table 7 Save and Export Options by Workflow Screen

Workflow screen	Optional actions	File type options	Export locations
Image preview	• Save a high-resolution image (HI Res image) • Export image information • Rename image	– .iBA – 16 Bit TIFF_Raw – 8 Bit Tiff – JPG	– USB – Network Drive
Contrast Versions	• Save a high-resolution image (HI Res image) • Export image information • Save a color image • Save all contrast versions • Rename image	– .iBA – 16 Bit TIFF_Raw – 8 Bit Tiff – JPG	– USB – Network Drive
Edit	• Save a high-resolution image (HI Res image)	– 8 Bit Tiff – JPG	– USB – Network Drive

Table 7 Save and Export Options by Workflow Screen (continued)

Workflow screen	Optional actions	File type options	Export locations
Gallery	1. Select the image to export. 2. Touch Export. • Rename image (one at a time). • Export image information	– .iBA – 16 Bit TIFF_Raw – 8 Bit Tiff	– USB – Network Drive

Additional information

When exporting images, select an appropriate file type for your application:

Table 8 Descriptions of File Types

File type	Description
16 Bit TIFF_Raw	High quality raw image, very large file size. Recommended for external image analysis and post-processing.
8 Bit Tiff	Medium-quality, large file suitable for standard analysis and documentation.
JPG	Medium-quality, small file suitable for presentations and reports.
.iBA	Proprietary file format for iBright™ Analysis Software. High-quality, very large file.
All Contrast Versions	Saves three grayscale contrast (Low, Medium, and High) for each image using the selected file format (JPG or TIFF). Note: Contrast Version are labeled Version A, B, C when the image was captured with a predefined Method using a manually set contrast stretch.
Hi Res image	Saves a high-resolution version of the image.
Color image	Saves a color image from the selected contrast version for publication purposes. The color image is exported as a 24-bit image.

Table 9 Descriptions of Export Locations

Locations	Description
USB drive	Selected image files can be saved to the USB drive. Subfolders can be created directly on the instrument to organize files.
Network drive	Saves image files to a connected network drive.



Band excision

The **Band Excision** feature allows users to identify, mark, and prepare specific gel bands for physical excision or downstream analysis.

IMPORTANT! Read the Radiation Safety section before performing band excision using UV or Blue light.

IMPORTANT! Do not use abrasive tools (e.g., razor blades, metal spatulas or metal tweezers) on the imaging trays, as they may scratch or damage the surface.

Excise bands from gels requiring UV illumination

1. From the **Home** screen, select **Band Excision** and follow the on-screen prompts
2. Wear appropriate personal protective equipment (PPE) including eye and skin protection.
3. Open the instrument drawer.
4. Place the safety guard and ensure that it is securely positioned.



WARNING! Bystanders without appropriate PPE must stay more than 1.5m (5ft) away. Failure to follow these precautions may result in serious eye and/or skin injury.



AVERTISSEMENT ! Les personnes ne portant pas d'équipement de protection individuelle (EPI) approprié doivent se tenir à une distance supérieure à 1,5 m (5 ft). Le non-respect de ces consignes peut entraîner des lésions graves aux yeux et/ou à la peau.

5. Confirm to turn the **UV light source** on.
6. Excise the bands as desired.
7. The light remains on for up to **05:00 minutes**.

Note: The light source turns off automatically when the countdown ends.

8. If required, use the light switch to manually turn the light source on or off.
9. Adjust the light output using the Intensity scale bar as needed.
 - Available intensity levels: 25%, 50%, 75%, 100%

10. Select **Back (<)** button to return to the Home screen.



WARNING! If the **Safety guard** is displaced, the light source automatically turns off, and the system returns to the Home screen.



AVERTISSEMENT ! Si le dispositif de protection est déplacé, la source lumineuse s'éteint automatiquement et le système revient à l'écran d'accueil.

Excise bands from gels requiring Blue illumination

1. From the **Home** screen, select **Band Excision** and follow the on-screen prompts
2. Wear appropriate PPE including eye and skin protection.
3. Open the instrument drawer.
4. Place the safety guard and ensure that it is securely positioned.



WARNING! Bystanders without appropriate PPE must stay more than 1.5m (5ft) away. Failure to follow these precautions may result in eye injury.



AVERTISSEMENT ! Les personnes ne portant pas d'équipement de protection individuelle (EPI) approprié doivent se tenir à une distance supérieure à 1,5 m (5 ft). Le non-respect de ces consignes peut entraîner des lésions oculaires.

5. Confirm to turn the **Blue light source** on .
6. Excise the band as desired.
7. The light remains on for up to **05:00 minutes**.

Note: The light source turns off automatically when the countdown ends.

8. If required, use the light switch to manually turn the light source on or off.
9. Adjust the light output using the Intensity scale bar as needed.
 - Available intensity levels: 25%, 50%, 75%, 100%
10. Select **Back (<)** button to return to the Home screen.

Note: If the **Safety guard** is displaced, the light source automatically turns off, and the system returns to the Home screen.

Excise bands from gels requiring white illumination

1. From the **Home** screen, select **Band Excision** and follow on-screen prompts
2. Wear appropriate PPE including eye and skin protection.
3. Open the instrument drawer.
4. Confirm to turn the **White light source** on or cancel to retrun to the home screen.
5. Excise the bands as desired.
6. The light remains on for up to **05:00 minutes**.
7. Use the light switch to manually turn the light source on or off.



Instrument maintenance

Follow the recommended maintenance schedule to prevent operational issues and ensure reliable results.

Administrator settings

To access **Settings**, from the **Home** screen, touch **Settings**.

The following categories are available: **Instrument**, **System**, and **Permissions**.

Table 10 Administrator Settings Description

Settings Category	Sub-Option	Description
Instrument	Network	Displays instrument details, including serial number, IP address, subnet mask, and firmware version. Allows execution of a manual self-test .
		• Configure network connectivity: – Wired: Connect using an Ethernet cable – Wireless: Connect using Wi-Fi (USB Wi-Fi Module, Cat. No. A26774 required) • Allows enabling or disabling Wi-Fi connectivity
	Time	• Set the time zone, date format, and time format for the region. • Allows synchronization of date and time when connected to the internet.
	Setting Control	• Enable or disable the on-screen keyboard. • Enable or disable network access. • Select system language. • Perform Flat-field correction (requires the White Tray with no sample present). • Perform a software update. • Enable automatic full-disk virus scan (infected files detected during scan will be automatically deleted)

Table 10 Administrator Settings Description (continued)

Settings Category	Sub-Option	Description
System	Function	• Enable or disable account creation. • Set automatic sign-out timer (default: 30 minutes). • Set sleep mode timer (default: 30 minutes). • Enable or disable access to the Gallery. • Define a gallery capacity threshold (for example 75%) and trigger a warning message when the gallery exceeds this limit.
	Module Parameters	• Set the default field of view (FOV) for image capture for each module (UV, Blue, and White).
	Export Options	• Configure default export options displayed during image saving, including image resolution, file format, export location, and DPI .
	Export Locations	• Add or delete network drive locations.
Permissions	Account Management	• Add or delete user accounts. • Edit user accounts, including resetting forgotten passwords.
	Method Management	• Create and manage predefined methods for each module. Note: Methods created by the administrator are visible to all users but editable only by the administrator. Note: Standard users can view method details but cannot modify them.
	Password Policy	• Define password composition requirements, including: – Uppercase letters – Lowercase letters – Numbers – Special characters • Configure: – Minimum password length – Password validity period (months) – Password attempt limit – Password history (retention of previous passwords)

Standard users settings

Access Settings from the "**Home**" screen.

Three options are available:

- Instrument
- System
- Permissions

Table 11 Standard Users Settings Description

Settings Category	Sub-Option	Description
Instrument	Network	Displays instrument details, including serial number, IP address, subnet mask, and firmware version. Allows execution of a manual self-test .
		• Configure network connectivity: – Wired: Connect using an Ethernet cable – Wireless: Connect using Wi-Fi (USB Wi-Fi Module, Cat. No. A26774 required) • Allows enabling or disabling Wi-Fi connectivity.
System	Export Locations	• Add or delete network drive locations.
Permissions	Account Management	• Edit user accounts, including changing passwords.

Run the instrument manual self-test

Instrument functions can be checked by a user to confirm proper hardware operating conditions.

1. From the Home screen Select Settings.
2. Select **Manual Self-Test**.
3. Confirm to Start Manual Self-Test.
4. During the self-test, the instruments check following:
 - a. Illumination modules
 - b. Motor operation
 - c. Camera initialization and shutter operation
 - d. Default system configurations

5. When the self-test is complete, a pop-up message displays one of the following statuses:
 - a. Self-test successful
 - b. Self-test failed

Note: If self test fails, contact tech support (thermofisher.com/support).

Clean the instrument

Repeated instrument use can cause spots and smudges on the sample trays, which can then decrease image quality. Clean the sample trays as needed.

Materials required

- Safety glasses
- Powder-free gloves
- Tissue, lint-free
- Deionized water
- Ethanol, 70% solution

Note: Avoid the use of detergents or abrasive cleaning agents.

1. Open the drawer to expose the sample tray.
2. Remove the sample tray.
3. Lightly spray the tray surface with deionized water or a 70% ethanol solution.
4. Wipe the surface with a lint-free tissue until sufficiently clean.
5. Clean the light box beneath the tray using 70% ethanol solution.
6. Close the drawer and operate the instrument as normal.



CAUTION! Do not spray liquids directly into the instrument or near electrical components. Apply cleaning solutions only to the sample tray and light box surfaces.



MISE EN GARDE ! Nettoyage et décontamination. Ne pulvérisez pas de liquides directement à l'intérieur de l'instrument ni à proximité des composants électriques. Appliquez les solutions de nettoyage uniquement sur le plateau d'échantillons et les surfaces du boîtier lumineux.



Instrument specifications

Instrument dimensions and specifications

Table 12 Dimensions

Specification	iGlow™ imager
Dimension (L x W x H)	38 × 45 × 41 cm (14.97 × 17.77 × 16.14 in)
Weight	22 kg (48.50lbs)
Instrument clearance (Back)	10 cm (3.94 in.)

Electrical requirements



CAUTION! For safety, the power outlet used for powering the instrument must be accessible at all times. In case of emergency, you must be able to immediately disconnect the main power supply to the instrument. Allow adequate space between the wall and the equipment so the power cord can be disconnected in case of emergency.



MISE EN GARDE ! Pour des raisons de sécurité, la prise de courant utilisée pour alimenter l'instrument doit être accessible à tout moment. En cas d'urgence, vous devez pouvoir déconnecter immédiatement l'alimentation principale de l'instrument. Prévoyez un espace suffisant entre le mur et l'équipement afin que le cordon d'alimentation puisse être débranché en cas d'urgence.

System electrical requirements:

Note: Do not plug this unit into a circuit shared with equipment that frequently power cycles (for example, refrigerators or freezers), as this may cause low-voltage brownout conditions and affect performance.

- Electric receptacle with grounding capability
- Maximum power dissipation: Approximately 168 W (not including computer and monitor)
- Mains AC line voltage tolerances must be within ± 10 percent of nominal voltage

Table 13 Electrical requirements

Device	Rated voltage	Circuit required	Rated frequency	Rated power	Static Power Consumption	Dynamic Power Consumption
iGlow™ Gel Documentation System (Ref: IG0001)	110–120 VAC	10 A	50/60 Hz	170 W	35 W	110 W
iGlow™ Gel Documentation System (Ref: IG0002)	200–240 VAC	10 A	50/60 Hz	170 W	35 W	110 W

Environmental requirements

Table 14 Environmental Requirements

Condition	Acceptable Range
Installation site	Indoor use only
Electromagnetic interference	Do not use this device in close proximity to sources of strong electromagnetic radiation (for example, unshielded intentional RF sources). Strong electromagnetic radiation may interfere with the proper operation of the device.
Altitude	Between sea level and 2000 m (6500 ft.) above sea level
Operating conditions	- Humidity: 15 – 80% relative humidity (noncondensing) - Temperature: 5 to 40°C (42 to 104°F) Note: For optimal performance, avoid rapid or extreme fluctuations in room temperature.
Storage and transport conditions	- Humidity: 20 – 80% relative humidity (noncondensing) - Temperature: –20 to 55°C (–4 to 131°F) - Atmospheric pressure: 75kPa to 106kPa
Thermal output	N/A
Vibration	Ensure that the instrument is not adjacent to strong vibration sources, such as a centrifuge, pump, or compressor. Excessive vibration will affect instrument performance.
Other conditions	Ensure the instrument is located away from any vents that could expel particulate material onto the instrument components. Avoid placing the instrument and computer adjacent to heaters, cooling ducts, or in direct sunlight.
Pollution degree	2

Network requirements

The Instrument is:

- Is factory-configured for IPv4 TCP/IP communication and includes an Ethernet adapter (100/1,000 Mbps) with an RJ45-type connector for integrating the device into a local area network (LAN)
- Can alternatively be configured for wireless networking (High Power USB WiFi Module required, sold separately as an optional accessory)

The instrument can be configured for *either* wired or wireless networking, not both.

Table 15 Required Materials

Network connection	Requirement
Wired	• Ethernet cable of sufficient length with RJ45 connectors • CAT5 cable for a 100 Mbps network
Wireless	High Power USB WiFi Module (Cat. No. A26774 , sold separately)

Instrument connections

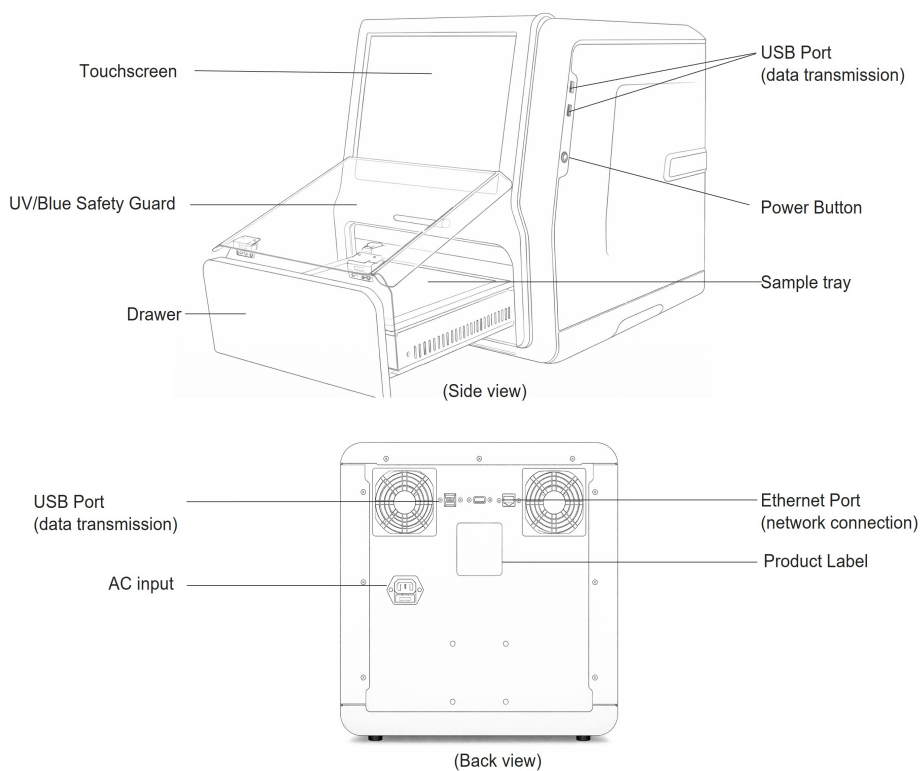


Figure 1 Instrument Connections



Safety



WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, and so on). To obtain SDSs, visit thermofisher.com/support.



AVERTISSEMENT ! SÉCURITÉ GÉNÉRALE. L'utilisation de ce produit d'une manière non conforme aux instructions figurant dans la documentation utilisateur peut entraîner des blessures corporelles ou endommager l'instrument ou l'appareil. Veillez à ce que toute personne utilisant ce produit soit formée aux bonnes pratiques de sécurité en laboratoire et ait pris connaissance des informations de sécurité fournies dans ce document.

- Avant toute utilisation d'un instrument ou d'un appareil, lisez attentivement et assurez-vous de bien comprendre les informations de sécurité figurant dans la documentation utilisateur fournie par le fabricant.
- Avant toute manipulation de produits chimiques, consultez et comprenez l'ensemble des fiches de données de sécurité (FDS) applicables et utilisez un équipement de protection individuelle approprié (gants, blouse, protection oculaire, etc.). Pour obtenir les FDS, consultez : thermofisher.com/support.

Symbols on this instrument

Symbols may be found on the instrument to warn against potential hazards or convey important safety information. In this document, the hazard symbol is used along with one of the following user attention words.

- **CAUTION!**—Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.
- **WARNING!**—Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.
- **DANGER!**—Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury.

Standard safety symbols

Symbol	English	Français
	Caution, risk of danger. Consult the manual for further safety information.	Attention, risque de danger. Consultez le manuel pour obtenir des informations de sécurité supplémentaires.
	Risk of electrical shock.	Risque de choc électrique.

Control and connection symbols

Symbols	Descriptions
	On (Power)
	Off (Power)
	Earth (ground) terminal
	Protective conductor terminal (main ground)
	Direct current
	Alternating current
	Both direct and alternating current

Conformity symbols

Conformity mark	Description
	Indicates conformity with safety requirements for Canada and U.S.A.
	Indicates conformity with China RoHS requirements.

(continued)

Conformity mark	Description
	Indicates conformity with Australian standards for electrical safety and electromagnetic compatibility.
	INDICATES CONFORMITY WITH THE WEEE DIRECTIVE 2012/19/EU.  CAUTION! To minimize negative environmental impact from disposal of electronic waste, do not dispose of electronic waste in unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provision and contact customer service for information about responsible disposal options.  MISE EN GARDE ! ÉLIMINATION DES DÉCHETS ÉLECTRONIQUES. Afin de réduire l'impact environnemental lié à l'élimination des déchets d'équipements électriques et électroniques, ne les jetez pas avec les déchets municipaux non triés. Respectez les réglementations locales en matière de gestion des déchets pour une élimination appropriée et contactez le service client pour obtenir des informations sur les solutions d'élimination responsables.
	Indicates conformity with European Union Requirements.

Safety information for instruments not manufactured by Thermo Fisher Scientific

Some of the accessories provided as part of the instrument system are not designed or built by Thermo Fisher Scientific. Consult the manufacturer's documentation for the information needed for the safe use of these products.

Instrument safety

General



CAUTION! Do not remove instrument protective covers. If you remove the protective instrument panels or disable interlock devices, you may be exposed to serious hazards including, but not limited to, severe electrical shock, laser exposure, crushing, or chemical exposure.

If covers are removed, do not use the instrument. Contact Technical Support.



MISE EN GARDE ! SÉCURITÉ DE L'INSTRUMENT. Ne retirez pas les capots de protection de l'instrument. Le retrait des panneaux de protection ou la neutralisation des dispositifs d'interverrouillage peut vous exposer à des dangers graves, notamment des chocs électriques sévères, une exposition au laser, des risques d'écrasement ou une exposition à des substances chimiques.

Si les capots ont été retirés, n'utilisez pas l'instrument. Contactez l'assistance technique.

Physical injury



CAUTION! Moving and Lifting Injury. The instrument is to be moved and positioned only by the personnel or vendor specified in the applicable site preparation guide. Improper lifting can cause painful and permanent back injury.

Things to consider before lifting or moving the instrument or accessories:

- Depending on the weight, moving or lifting may require two or more persons.
- If you decide to lift or move the instrument after it has been installed, do not attempt to do so without the assistance of others, the use of appropriate moving equipment, and proper lifting techniques.
- Ensure you have a secure, comfortable grip on the instrument or accessory.
- Make sure that the path from where the object is to where it is being moved is clear of obstructions.
- Do not lift an object and twist your torso at the same time. Keep your spine in a good neutral position while lifting with your legs.
- Participants should coordinate lift and move intentions with each other before lifting and carrying.
- For smaller packages, rather than lifting the object from the packing box, carefully tilt the box on its side and hold it stationary while someone else slides the contents out of the box.



MISE EN GARDE ! RISQUE LIÉ AU DÉPLACEMENT ET AU LEVAGE. L'instrument doit être déplacé et installé uniquement par le personnel ou le prestataire spécifié dans le guide de préparation du site applicable. Un levage inapproprié peut entraîner des blessures dorsales douloureuses et permanentes.

Points à prendre en compte avant de déplacer ou de soulever l'instrument ou ses accessoires:

- Selon le poids, le déplacement ou le levage peut nécessiter deux personnes ou plus.
- Si vous devez déplacer ou soulever l'instrument après son installation, ne le faites pas sans assistance, sans équipement de manutention approprié et sans appliquer des techniques de levage adaptées.
- Assurez-vous d'avoir une prise sûre et confortable sur l'instrument ou l'accessoire.
- Vérifiez que le trajet à parcourir est dégagé de tout obstacle.
- Ne soulevez pas une charge en effectuant une rotation du torse. Maintenez la colonne vertébrale en position neutre et soulevez en utilisant les jambes.
- Les personnes impliquées doivent se coordonner avant de soulever et de transporter la charge.
- Pour les colis de petite taille, plutôt que de soulever l'objet hors de son emballage, inclinez soigneusement la boîte sur le côté, maintenez-la en position pendant qu'une autre personne en fait glisser le contenu.

Electrical safety



WARNING! Fuse Installation. Before installing the instrument, verify that the fuses are properly installed and the fuse voltage matches the supply voltage. Replace fuses only with the type and rating specified for the unit. Improper fuses can damage the instrument wiring system and cause a fire.



AVERTISSEMENT ! Installation des fusibles. Avant d'installer l'instrument, vérifier que les fusibles sont correctement insérés et que leur tension correspond à celle fournie par le circuit d'alimentation. Ne remplacer les fusibles que par des modèles du type et de la puissance spécifiés pour l'appareil. L'utilisation de fusibles inadaptés peut endommager le circuit électrique de l'instrument et provoquer un incendie.



WARNING! Ensure appropriate electrical supply. For safe operation of the instrument:

- Plug the system into a properly grounded receptacle with adequate current capacity.
- Ensure the electrical supply is of suitable voltage.
- Never operate the instrument with the ground disconnected. Grounding continuity is required for safe operation of the instrument.



AVERTISSEMENT ! Veiller à utiliser une alimentation électrique appropriée. Pour garantir le fonctionnement de l'instrument en toute sécurité :

- Brancher le système sur une prise électrique correctement mise à la terre et de puissance adéquate.
- S'assurer que la tension électrique est convenable.
- Ne jamais utiliser l'instrument alors que le dispositif de mise à la terre est déconnecté. La continuité de la mise à la terre est impérative pour le fonctionnement de l'instrument en toute sécurité.



WARNING! Power Supply Line Cords. Use properly configured and approved line cords for the power supply in your facility. If the line cord is damaged, contact Technical Support.



AVERTISSEMENT ! Cordons d'alimentation électrique. Utiliser des cordons d'alimentation adaptés et approuvés pour raccorder l'instrument au circuit électrique du site.



WARNING! Disconnecting Power. To fully disconnect power either detach or unplug the power cord, positioning the instrument such that the power cord is accessible.



AVERTISSEMENT ! Déconnecter l'alimentation. Pour déconnecter entièrement l'alimentation, détacher ou débrancher le cordon d'alimentation. Placer l'instrument de manière à ce que le cordon d'alimentation soit accessible.

UV safety



WARNING! RISK GROUP 2. CAUTION UV emitted from this product. Do not stare at the lamp. May be harmful to eyes and skin.



AVERTISSEMENT ! GROUPE DE RISQUE 2. Attention aux rayons UV émis par ce produit. Ne regardez pas directement vers la lampe de fonctionnement. Il peut y avoir des dommages aux yeux et à la peau.

Cleaning and decontamination



CAUTION! Cleaning and Decontamination. Use only the cleaning and decontamination methods that are specified in the manufacturer user documentation. It is the responsibility of the operator (or other responsible person) to ensure that the following requirements are met:

- No decontamination or cleaning agents are used that can react with parts of the equipment or with material that is contained in the equipment. Use of such agents could cause a HAZARD condition.
- The instrument is properly decontaminated a) if hazardous material is spilled onto or into the equipment, and/or b) before the instrument is serviced at your facility or is sent for repair, maintenance, trade-in, disposal, or termination of a loan. Request decontamination forms from customer service.
- Before using any cleaning or decontamination methods (except methods that are recommended by the manufacturer), confirm with the manufacturer that the proposed method will not damage the equipment.



MISE EN GARDE ! Nettoyage et décontamination. Utiliser uniquement les méthodes de nettoyage et de décontamination indiquées dans la documentation du fabricant destinée aux utilisateurs. L'opérateur (ou toute autre personne responsable) est tenu d'assurer le respect des exigences suivantes:

- Ne pas utiliser d'agents de nettoyage ou de décontamination susceptibles de réagir avec certaines parties de l'appareil ou avec les matières qu'il contient et de constituer, de ce fait, un DANGER.
- L'instrument doit être correctement décontaminé a) si des substances dangereuses sont renversées sur ou à l'intérieur de l'équipement, et/ou b) avant de le faire réviser sur site ou de l'envoyer à des fins de réparation, de maintenance, de revente, d'élimination ou à l'expiration d'une période de prêt (des informations sur les formes de décontamination peuvent être demandées auprès du Service clientèle).
- Avant d'utiliser une méthode de nettoyage ou de décontamination (autre que celles recommandées par le fabricant), les utilisateurs doivent vérifier auprès de celui-ci qu'elle ne risque pas d'endommager l'appareil.

Instrument component and accessory disposal

To minimize negative environmental impact from disposal of electronic waste, do not dispose of electronic waste in unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provision and contact customer service for information about responsible disposal options.

Safety and electromagnetic compatibility (EMC) standards

The instrument design and manufacture complies with the following standards and requirements for safety and electromagnetic compatibility.

Safety and EMC standards

Reference	Description
EU Directive 2014/35/EU	European Union “Low Voltage Directive”
IEC 61010-1 EN 61010-1 UL 61010-1 CAN/CSA C22.2 No. 61010-1	<i>Safety requirements for electrical equipment for measurement, control, and laboratory use – Part 1: General requirements</i>
IEC 62471 EN 62471	<i>Photobiological safety of lamps and lamp systems</i>

EMC standards

Reference	Description
EU Directive 2014/30/EU	European Union “EMC Directive”
EN 61326-1 IEC 61326-1	<i>Electrical Equipment for Measurement, Control and Laboratory Use – EMC Requirements – Part 1: General Requirements</i>
AS/NZS CISPR 11	<i>Limits and Methods of Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radiofrequency Equipment</i>

(continued)

Reference	Description
FCC Part 15 Subpart B (47 CFR)	<i>U.S. Standard Radio Frequency Devices</i> This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense. This equipment has been designed and tested to CISPR 11 Class A. In a domestic environment it may cause radio interference, in which case, you may need to take measures to mitigate the interference. Do not use this device in close proximity to sources of strong electromagnetic radiation (e.g. unshielded intentional RF sources), as these can interfere with the proper operation.
ICES-003 Issue 7	Interference-Causing Equipment Standard (ICES) for Information Technology Equipment (including Digital Apparatus). CAN ICES-003(A) / NMB-003(A)

Environmental design standards

Reference	Description
Directive 2012/19/EU	European Union “WEEE Directive” — Waste electrical and electronic equipment
Directive 2011/65/EU & Commission Delegated Directive 2015/863	European Union “RoHS Directive” — Restriction of hazardous substances in electrical and electronic equipment
SJ/T 11364-2024	“China RoHS” Standard — Marking for the Restricted Use of Hazardous Substances in Electronic and Electrical Products For instrument specific certificates, visit our customer resource page at www.thermofisher.com/us/en/home/technical-resources/rohs-certificates.html.

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the "Documentation and Support" section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with sufficient ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if needed) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.



AVERTISSEMENT ! PRÉCAUTIONS GÉNÉRALES EN CAS DE MANIPULATION DE PRODUITS CHIMIQUES.

Pour minimiser les risques, veiller à ce que le personnel du laboratoire lise attentivement et mette en œuvre les consignes de sécurité générales relatives à l'utilisation et au stockage des produits chimiques et à la gestion des déchets qui en découlent, décrites ci-dessous. Consulter également la FDS appropriée pour connaître les précautions et instructions particulières à respecter :

- Lire et comprendre les fiches de données de sécurité (FDS) fournies par le fabricant avant de stocker, de manipuler ou d'utiliser les matériaux dangereux ou les produits chimiques. Pour obtenir les FDS, se reporter à la section « Documentation et support » du présent document.
- Limiter les contacts avec les produits chimiques. Porter des équipements de protection appropriés lors de la manipulation des produits chimiques (par exemple : lunettes de sûreté, gants ou vêtements de protection).
- Limiter l'inhalation des produits chimiques. Ne pas laisser les récipients de produits chimiques ouverts. Ils ne doivent être utilisés qu'avec une ventilation adéquate (par exemple, sorbonne).
- Vérifier régulièrement l'absence de fuite ou d'écoulement des produits chimiques. En cas de fuite ou d'écoulement d'un produit, respecter les directives de nettoyage du fabricant recommandées dans la FDS.
- Manipuler les déchets chimiques dans une sorbonne.

- Veiller à utiliser des récipients à déchets primaire et secondaire. (Le récipient primaire contient les déchets immédiats, le récipient secondaire contient les fuites et les écoulements du récipient primaire. Les deux récipients doivent être compatibles avec les matériaux mis au rebut et conformes aux exigences locales, nationales et communautaires en matière de confinement des récipients.)
- Une fois le récipient à déchets vidé, il doit être refermé hermétiquement avec le couvercle fourni.
- Caractériser (par une analyse si nécessaire) les déchets générés par les applications, les réactifs et les substrats particuliers utilisés dans le laboratoire.
- Vérifier que les déchets sont convenablement stockés, transférés, transportés et éliminés en respectant toutes les réglementations locales, nationales et/ou communautaires en vigueur.
- **IMPORTANT !** Les matériaux représentant un danger biologique ou radioactif exigent parfois une manipulation spéciale, et des limitations peuvent s'appliquer à leur élimination.



WARNING! HAZARDOUS WASTE (from instruments). Waste produced by the instrument is potentially hazardous. Follow the guidelines noted in the preceding General Chemical Handling warning.



AVERTISSEMENT ! DÉCHETS DANGEREUX (PROVENANT DE L'INSTRUMENT). Les déchets générés par l'instrument peuvent présenter un caractère dangereux. Respectez les consignes indiquées dans l'avertissement relatif à la manipulation générale des produits chimiques ci-dessus.

Biological hazard safety



WARNING! Potential Biohazard. Depending on the samples used on this instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards.



AVERTISSEMENT ! RISQUE BIOLOGIQUE POTENTIEL. Selon la nature des échantillons utilisés avec cet instrument, ses surfaces peuvent être considérées comme présentant un risque biologique. Appliquez des méthodes de décontamination appropriées lors de toute manipulation de matériaux biologiquement dangereux.



Documentation and support

Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.

